

PRODUCT INFORMATION



Mitragynine Pseudoindoxyl

Item No. 20033

Formal Name: ($\alpha E,6'S,7'S,8'aS$)-6'-ethyl-1,2',3,3',6',7',8',8'a-octahydro-4-methoxy- α -(methoxymethylene)-3-oxo-spiro[2H-indole-2,1'(5'H)-indolizine]-7'-acetic acid, methyl ester

MF: $C_{23}H_{30}N_2O_5$

FW: 414.5

Purity: $\geq 98\%$

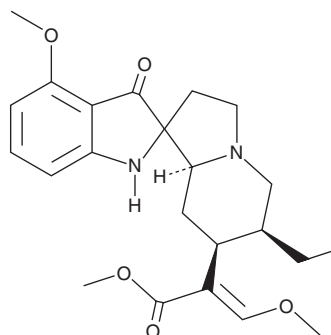
Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 2 years

Special Conditions: Light-sensitive

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.



Description

Mitragynine pseudoindoxyl (Item No. 20033) is an analytical reference standard categorized as an opioid.^{1,2} Mitragynine pseudoindoxyl has analgesic activity in mice without inducing the development of opioid-induced hyperalgesia.² It induces physical dependence in mice but reduces signs of withdrawal in morphine-dependent mice administered naloxone. This product is intended for research and forensic applications.

References

- Váradí, A., Marrone, G.F., Palmer, T.C., *et al.* Mitragynine/corynantheidine pseudoindoxyls as opioid analgesics with mu agonism and delta antagonism, which do not recruit β -arrestin-2. **59(18)**, 8381-8397 (2016).
- Wilson, L.L., Chakraborty, S., Eans, S.O., *et al.* Kratom alkaloids, natural and semi-synthetic, show less physical dependence and ameliorate opioid withdrawal. *Cell. Mol. Neurobiol.* **41(5)**, 1131-1143 (2021).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/28/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM